

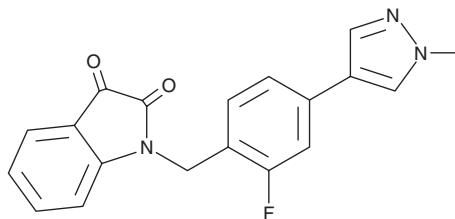
PRODUCT INFORMATION



ML-137

Item No. 40918

CAS Registry No.: 1222834-53-8
Formal Name: 1-[[2-fluoro-4-(1-methyl-1H-pyrazol-4-yl)phenyl]methyl]-1H-indole-2,3-dione
Synonym: VU0366369
MF: C₁₉H₁₄FN₃O₂
FW: 335.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ML-137 is supplied as a solid. A stock solution may be made by dissolving the ML-137 in the solvent of choice, which should be purged with an inert gas. ML-137 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

ML-137 is a positive allosteric modulator (PAM) of M₁ muscarinic acetylcholine receptors (mAChRs).¹ It is selective for M₁ over M₂-M₅ mAChRs at 30 μM. ML-137 potentiates ACh-induced activation of the M₁ muscarinic receptor (EC₅₀ = 0.83 μM).

Reference

1. Bridges, T.M., Phillip Kennedy, J., Noetzel, M.J., et al. Chemical lead optimization of a pan Gq mAChR M1, M3, M5 positive allosteric modulator (PAM) lead. Part II: Development of a potent and highly selective M1 PAM. *Bioorg. Med. Chem. Lett.* **20(6)**, 1972-1975 (2010).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/03/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM